

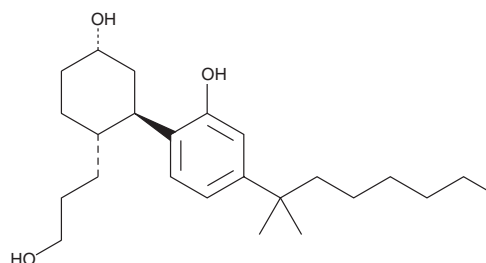
PRODUCT INFORMATION



(±)5-*epi* CP 55,940

Item No. 13803

CAS Registry No.: 2365471-91-4
Formal Name: *rel*-5-(1,1-dimethylheptyl)-2-
[(1*R*,2*R*,5*S*)-5-hydroxy-2-(3-
hydroxypropyl)cyclohexyl]-phenol
MF: C₂₄H₄₀O₃
FW: 376.6
Purity: ≥97%
UV/Vis.: λ_{max}: 219, 277 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

(±)5-*epi* CP 55,940 (Item No. 13803) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. (±)5-*epi* CP 55,940 is a potential byproduct in the synthesis of (±)-CP 55,940 (Item No. 13241) based on the published synthesis of (-)-CP 47,497 (Item No. 10917).¹⁻³ This product is intended for research and forensic applications.

References

1. Melvin, L.S., Johnson, M.R., Harbert, C.A., *et al.* A cannabinoid derived prototypical analgesic. *J. Med. Chem.* **27**(1), 67-71 (1984).
2. Johnson, M.R. and Lawrence S.M. Jr., J. 2-Hydroxy-4-(substituted) phenyl cycloalkanes and derivatives. *Pfizer, Inc.* **US4371720A** (1983).
3. Hengst, J.A., Nduwumwami, A.J., Raup-Konsavage, W.M., *et al.* Inhibition of sphingosine kinase activity enhances immunogenic cell surface exposure of calreticulin induced by the synthetic cannabinoid 5-*epi*-CP-55,940. *Cannabis Cannabinoid Res.* **7**(5), 637-647 (2022).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/08/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM